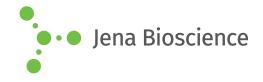
# **DATA SHEET**





#### Rsal

Isoschizomers: Afal, Neoschizomers: Csp6l JBSpeed Restriction Enzyme

Cat. No.	Amount
EN-129S	1.000 Units
EN-129L	5 x 1000 Units

5'	G	Т	<b>1</b>	Α	С	3'
3'	С	Α	<b>↑</b>	Т	G	5'

**Unit Definition:** One unit is the amount of enzyme required to completely digest 1  $\mu g$  of Lambda DNA (113 sites) in 1 hour in a total reaction volume of 50  $\mu$ l. Enzyme activity was determined in the recommended reaction buffer.

For general laboratory use.

**Shipping:** shipped on gel packs **Storage Conditions:** store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Form: liquid (Supplied in 10 mM Tris-HCl pH 7.4, 50 mM KCl, 0.1 mM

EDTA, 200 µg/ml BSA and 50 % [v/v] glycerol)

Concentration: 10 units/µl

Source: Rhodopseudomonas sphaeroides

Supplied with: 10x Universal Buffer (UB)

#### Recommended 50 µl assay

5 μl	10x Universal Buffer (UB)		
1 μg	pure DNA <sup>1</sup> or PCR product <sup>2</sup>		
10 units	enzyme		
fill up to 50 μl	PCR grade water		

<sup>&</sup>lt;sup>1</sup> Supercoiled or high molecular weight DNA (e.g. plant genomic DNA) may require longer incubation time or higher amount of enzyme.

#### **Protocol:**

- The enzyme should not exceed 10 % of total reaction volume.
- Add enzyme as last component. Mix components well before adding enzyme. After enzyme addition, mix gently by pipetting. Do not vortex.
- Incubate 5 to 10 min. at 37 °C.
- · Stop reaction by alternatively:
  - Addition of 2.1  $\mu$ l EDTA pH 8.0 [0.5 M], final 20 mM
  - Heat Inactivation (20 min. at 65 °C)
  - Spin Column DNA Purification (e.g. PCR Purification Kit, Cat.-No. PP-201S/L)
  - Gel Electrophoresis and Single Band Excision (e.g. Agarose Gel Extraction Kit, Cat.-No. PP-202S/L)
  - Phenol-Chloroform Extraction or Ethanol Precipitation.

#### **Double Digestion - Buffer Compatibility:**

B1 - 75-100 % Relative Activity

B2 - 100 % Relative Activity

B3 - 50 % Relative Activity

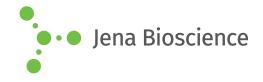
B4 - <10 % Relative Activity B5 - <10 % Relative Activity

1x UB - 100 % Relative Activity (recommended)

Please note that the optimum digestion condition for this enzyme is 1x UB. Within the Universal Buffer (UB) system, the most majority of our enzymes display 100% Relative Activity in 1x UB and only few either in 0.5x or 2x UB. If optimum condition for second enzyme is different than the recommended for the first enzyme, we suggest carrying out first the restriction at the higher recommended concentration of UB and dilute the reaction volume to the adequate UB concentration for further proceeding with the second restriction.

 $<sup>^{2}</sup>$  Some enzymes may require additional DNA bases flanking the restriction site for complete digestion.

# DATA SHEET





## **Rsal**

Isoschizomers: Afal, Neoschizomers: Csp6l JBSpeed Restriction Enzyme

#### **Reaction Enzymes Buffer Guide:**

Redection Enzymes burier dulue.						
Buffer 1	10 × B1	100 mM 100 mM 1000 μg/ml	Tris-HCl (pH 7.9, 25°C) MgCl <sub>2</sub> BSA			
Buffer 2	10 × B2	100 mM 100 mM 500 mM 1000 µg/ml	Tris-HCl (pH 7.9, 25°C) MgCl <sub>2</sub> NaCl BSA			
Buffer 3	10 × B3	500 mM 100 mM 1000 mM 1000 µg/ml	Tris-HCl (pH 7.9, 25°C) MgCl <sub>2</sub> NaCl BSA			
Buffer 4	10 × B4	100 mM 100 mM 1500 mM 1000 µg/ml	Tris-HCl (pH 7.9, 25°C) MgCl <sub>2</sub> NaCl BSA			
Buffer 5	10 × B5	200 mM 100 mM 500 mM 1000 µg/ml	Tris-acetate (pH 7.9, 25°C) Mg-acetate K-acetate BSA			

# Reaction Buffer Compatibility:

Our restriction enzymes are fully compatible to restrictases and buffer systems from other manufacturers and can be used along in double digestions. To obtain best results, consult the corresponding manuals of all involved products.

#### **Ligation and recutting:**

After 10-fold overdigestion with Rsal, >95% of the DNA fragments can be ligated and recut with this enzyme.

#### **DNA Methylation:**

No Inhibition: dam, dcm Inhibition (Blocked by some combinations of overlapping): CpG

### **Quality Control:**

All preparations are assayed for contaminating endonuclease, 3'-exonuclease, 5' exonuclease/ 5' phosphatase, as well as nonspecific single- and doublestranded DNase activities.

#### **Related Products:**

Universal Restriction Buffer, #EN-300